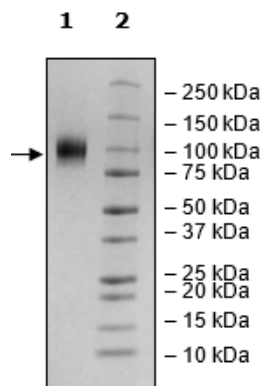


Product Information

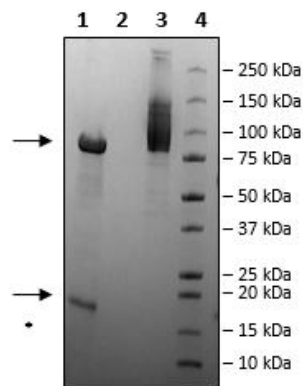
Description:	Recombinant human CD4, encompassing amino acids 26-396, corresponding to the extracellular domain. This construct includes an Fc-domain from IgG1 fused at the C-terminus, followed by an Avi-tag™. This protein was affinity purified.
Background:	CD4 is a cell surface glycoprotein found on partially defined functional T cell subsets, including helper T cells and T-regulatory cells, peripheral monocytes and other APCs (antigen-presenting cells). The ectodomain of CD4 binds to membrane-proximal domains of MHC (major histocompatibility complex) class II molecules, while its cytoplasmic domains interact with the protein tyrosine kinase p56lck (lck) through a shared cysteine-containing motif. CD4 ⁺ T cells, or helper T cells, are a type of lymphocyte that helps coordinate the immune response against infection and disease by activating cells of the innate immune system, B-lymphocytes and cytotoxic T cells. CD4 ⁺ T cells are activated by interaction between the TCR (T- cell receptor) and its cognate peptide presented on MHC II molecules, and CD4 is a critical component of the T cell receptor complex that recognizes antigenic peptides presented by MHC II molecules, increasing its stability. CD4 is typically used as a T cell marker that allows to characterize the populations of T cells present in a sample.
Species:	Human
Construct:	CD4 (26-396-Fc(IgG1)-Avi)-(Biotin)
Concentration:	1.24 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	70 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_000616.5
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful for SDS-PAGE and avidin-pull down assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown



1. Beads
2. Flow thru
3. Control
4. Standards

* Avidin from beads.